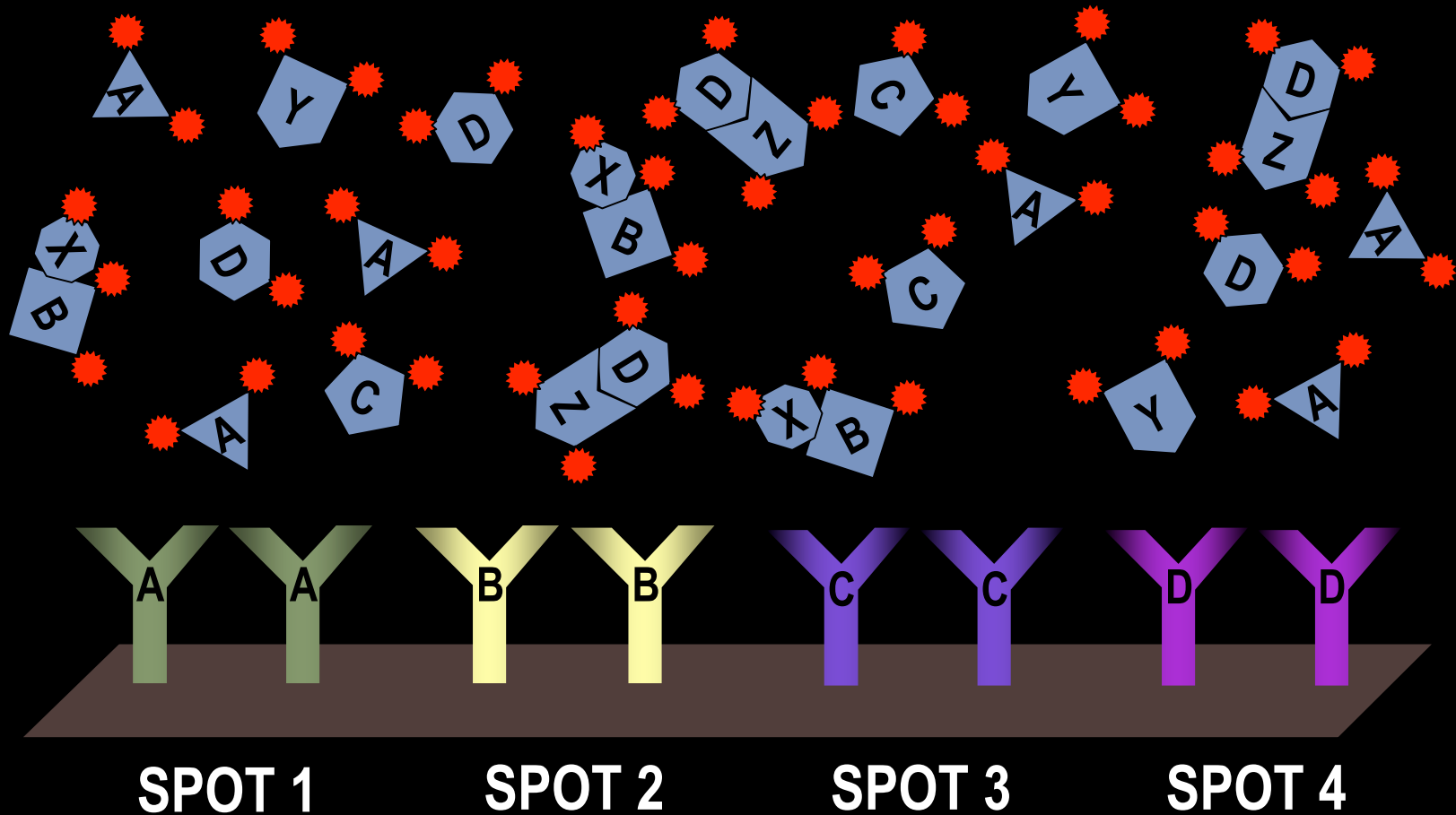


# Kinex™ Antibody Microarrays

Step 3



KINEXUS



[www.kinexus.ca](http://www.kinexus.ca)